



BIO-RAD

Declaration of Conformity

Bio-Rad Laboratories, Inc., 1000 Alfred Nobel Drive, Hercules, California, 94547, U.S.A., declares that the product

175 Tube Cell

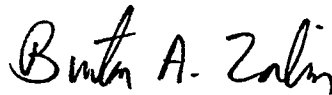
to which this declaration relates, is in conformity to the following standards or normative documents

EN55011:1991
EN50082-2:1994
EN61000-4-2:1995
ENV50141:1994
ENV50140:1994
EN61000-4-4:1995
EN61000-4-8:1995
EN61010-1:1993

following the provisions of the 89/336/EEC & 73/23/EEC Directive.

The product is imported into the EU by Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead (London area), Hertfordshire HP2 7TD England.

29 November, 1995
date of issue



Burton Zabin
Vice President

DC1651980 Rev A